

## Bio-Scale DEAE20 Anion Exchange Column

(15 mm x 118 mm)    Catalog Number 751-0037

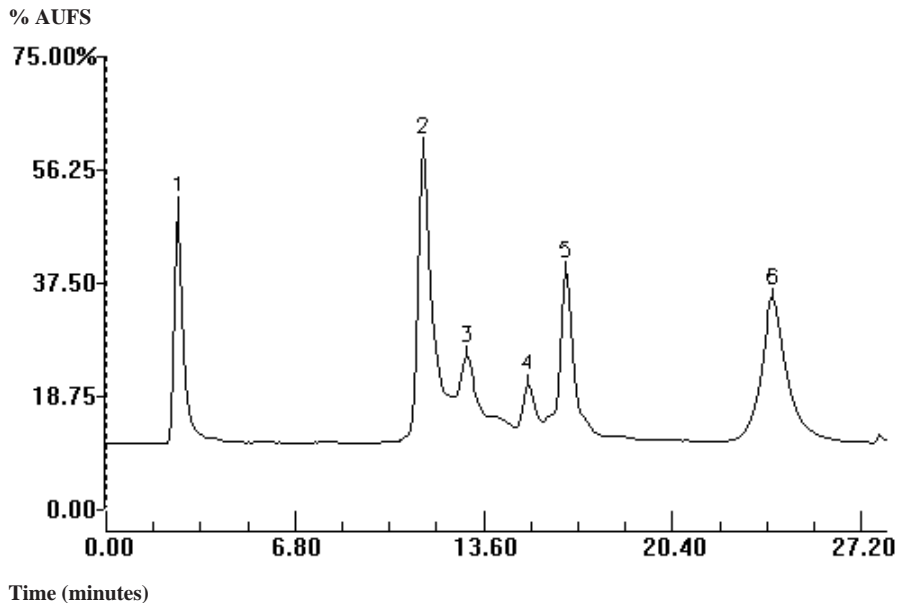
Serial Number: \_\_\_\_\_  
Batch Number: \_\_\_\_\_  
Resin Batch #: \_\_\_\_\_  
Date Tested: \_\_\_\_\_  
  
Checked By: \_\_\_\_\_

### Column Specifications

Efficiency (N):	$\geq 12,000$ / meter
Asymmetry (b/a):	0.85 - 1.50
Column Pressure (4.59 ml/min):	< 160 psi

### Typical Chromatogram\*

<b>Buffer A:</b> 20 mM Tris, pH 7.6	<b>Sample:</b> Protein Standard for Anion Exchange Chromatography (Catalog Number 125-0561)
<b>Buffer B:</b> 20 mM Tris + 1.0 M Sodium Chloride pH 7.6	
<b>Gradient:</b> 0% B in 6 minutes 0 - 50%B in 20 minutes	<b>Inj. Vol:</b> 1000 $\mu$ l (9000 $\mu$ g Total Protein)
<b>Flow Rate:</b> 4.59 ml/min	<b>Peaks:</b> 1. Myoglobin 2. Conalbumin - A 3. Conalbumin - B 4. Ovalbumin 5. Soybean Trypsin Inhibitor
<b>Detection:</b> UV @ 280 nm, 0.16 AUFS	



\* The above chromatogram and test conditions are provided to help resolve any technical problems.